

LANCE® Ultra

Eu-W1024 labeled Anti-phospho-Crosstide (GSK-3 α Ser21)**Antibody**

Product number: TRF0202-D **Lot Number:** 3251202

Product Format: TRF0202-D: 10 μ g
TRF0202-M: 100 μ g

Manufacturing date: 9/26/2023 **Document version:** 1

Product Information

Antibody: Europium-labeled mouse monoclonal antibody recognizing phospho-Ser21 in peptides derived from glycogen synthase kinase-3 α (GSK-3 α)

Storage Buffer: 50 mM Tris-HCl (pH 7.4), 0.9% NaCl with 0.05% sodium azide as preservative and 0.1 % BSA

Molecular Weight: 160 000

Stability: This product is stable for at least 24 months from the manufacturing date when stored in its original packaging and at the recommended storage conditions.

Storage Conditions: Store at 4°C.

Quality Control

The QC release specifications are based on spectrophotometric analysis of the labeled antibody. We certify that results meet our quality release criteria.

Labeling Ratio: 5.23
Concentration: 100 μ g/mL (0.625 μ M)

Recommended Assay Conditions

MSK1: ATP titration

SUGGESTED METHOD:

(Specific applications might require optimization)

Reagent Preparation

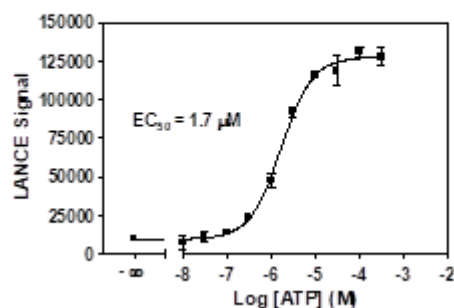
- Prepare 1X Kinase Assay Buffer: 50 mM HEPES pH 7.5, 1 mM EGTA, 10 mM MgCl₂, 3 mM MnCl₂, 2 mM DTT and 0.01% Tween-20.
- Prepare a 2X MSK1 solution: dilute the enzyme to a concentration of 4 nM in Kinase Assay Buffer. Keep on ice.
- Prepare a 4X mix of *ULight*[™]-Crosstide: dilute *ULight*[™]-Crosstide to a concentration of 200 nM in Kinase Assay Buffer.
- Prepare a 4X mix of ATP: dilute ATP to concentrations ranging from 40 nM to 4 mM (serial half-log dilutions) in Kinase Assay Buffer. Keep on ice.
- Prepare 1X Detection Buffer: dilute 0.5 mL of 10X Detection Buffer with 4.5 mL of H₂O.
- Prepare a 4X Stop Solution: prepare a 40 mM EDTA solution in 1X Detection Buffer.
- Prepare a 4X Detection Mix: dilute the Eu-anti-phospho-Crosstide antibody to a concentration of 8 nM in 1X Detection Buffer.

Protocol

- Pipet 5 µL of 2X MSK1 solution into a 384-well white OptiPlate[™]-384 (2 nM final concentration).
- Add 2.5 µL of 4X *ULight*[™]-Crosstide (50 nM final concentration).
- Add 2.5 µL of 4X ATP mix (1 mM to 10 nM final concentrations).
- Cover plate with TopSeal-A and incubate 60 min at 23°C.
- Add 5 µL of 4X Stop Solution and incubate 5 min at 23°C.
- Add 5 µL of Detection Mix (2 nM Eu-anti-phospho-Crosstide antibody final concentration) and mix.
- Cover plate with TopSeal-A and incubate 60 min at 23°C.
- Remove TopSeal-A and read in TR-FRET mode at 665 nm (excitation at 320 or 340 nm).

Typical Product Data

MSK1 kinase assay using ULight™-Crosstide and E u-anti-phospho-Crosstide antibody obtained using the EnVision® Multilabel Reader:



Suggested Materials

	Supplier
• Substrate: ULight™- Crosstide	Revvity Inc
• Antibody: Eu- anti-phospho-Crosstide (GSK-3α Ser21)	Revvity Inc
• Kinase: MSK1, active	Millipore/Upstate
• Detection Buffer: LANCE® Detection Buffer, 10X	Revvity Inc
• Plate: OptiPlate™-384, white	Revvity Inc
• TopSeal™-A	Revvity Inc

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

TRF0202-R Rev01

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.